



## Anti-CRTH2 (IN)

*(Chemoattractant receptor-homologous molecule expressed on Th2 cells, GPR44)*

**CATALOG NO.: 54647**

### BACKGROUND:

The chemoattractant receptor-homologous molecule expressed on Th2 cells (CRTH2) is a recently identified receptor for the prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) (1,2) in addition to the classic prostaglandin D receptor. CRTH2 is expressed on Th2 cells and eosinophils and mediates chemotaxis of these cells to PGD<sub>2</sub> and is thus thought to be a key receptor mediating eosinophil and Th2 recruitment during allergic responses. However, CRTH2-null mice showed enhanced eosinophil recruitment into the lung consistent with observations that the CRTH2-null mice produced significantly higher amounts of interleukin-5 (IL-5) and IL-3 (3). This suggests that CRTH2 plays a nonredundant role in restricting eosinophilia and allergic response in vivo. At least two different isoforms of CRTH2 are known to exist.

### SOURCE & REACTIVITY:

Rabbit polyclonal anti-CRTH2 was raised against a 15 amino acid peptide from near the center of human CRTH2 (GenBank accession no. NP\_034092). Anti-CRTH2 is human, mouse, and rat reactive.

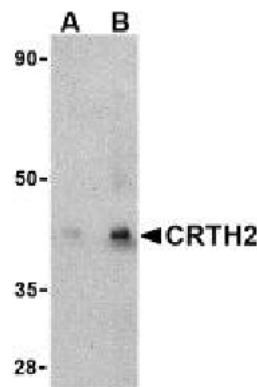
### APPLICATION:

The following concentration ranges are recommended starting points for this product.

### Western Blot:

2.0 µg/ml

Positive Control: Human heart tissue lysate



Western blot analysis of CRTH2 in human heart tissue lysate with anti-CRTH2 at (A) 1 and (B) 2 µg/ml.

*This product is for in vitro research purposes only.*

### RELATED PRODUCTS:

Anti-CRTH2 (NT), Catalog No. **54646**

### STORAGE:

The antibody is supplied as purified IgG, 50 µg in 250 µl of 1X PBS containing 0.02% sodium azide. Store at 4 °C for up to one year. Avoid repeated freezing and thawing.

### REFERENCES:

1. Nagata K, et al (1999) *FEBS Lett.* 459:195-9.
2. Shichijo M, et al (2003) *J. Pharmacol. Exp. Ther.* 307:518-25.
3. Chevalier E, et al (2005) *J. Immunol.* 2056-60.